

Product Number:	100860
Lot Number:	502293
Expiration:	3-Sep-2020
Platform(s):	HD-1/ HD-X



Data below represents results generated on the Simoa™ HD-1 Analyzer.

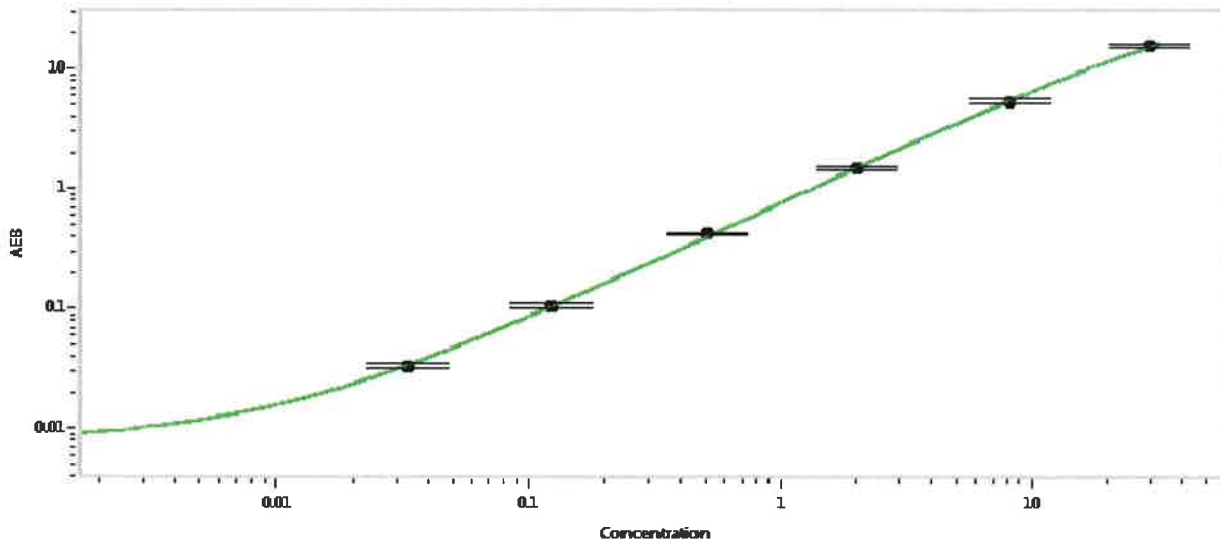
Component	Part Number	Lot Number
Bead Reagent	100865	926805
Detector Reagent	100862	004504
SBG Reagent	100868	924612
Sample Diluent	100870	924610
Calibrators	101111	926101
Controls	100875 & 100876	926102
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	926102	5.48	4.42-6.63
Control 2	926102	0.080	0.069-0.103
Panel 1	908501	1.09	0.907-1.99
Panel 2	908502	6.37	5.23-12.7
Panel 3	908503	20.4	15.5-47.6

* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.033
C	0.124
D	0.512
E	2.03
F	8.20
G	30.1

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>p stadler</i>	20 Mar 2020
	Name	Title	Signature/Date	